

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-Plex® and V-Plex Plus Human IL-10 Kit
Kit Catalog Numbers: K151QU-Series
Kit Lot Number: K0081078
Expiration Date: 31 Jan 2020

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-10	0.298	80% - 120%	< 20%	233	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-10	Calibrator	359	pg/mL
		Proinflammatory Panel 1 (human) Control 1	125	pg/mL
		Proinflammatory Panel 1 (human) Control 2	24.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.47	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
359	186,000	1.0
89.8	48,300	1.5
22.4	11,200	2.0
5.61	2,770	2.7
1.40	764	1.3
0.351	243	1.7
0.0876	117	4.2
0.00	75	4.8

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-10 Plate	Z0046903	2-8°C	28 Feb 2021
Anti-hu IL-10 Antibody	D0081024	2-8°C	31 Jan 2020
Proinflammatory Panel 1 (human) Calibrator Blend	A0080174	2-8°C	31 Jul 2023
Diluent 2	M0020099	≤ -10°C	31 Mar 2020
Diluent 3	M0030084	≤ -10°C	31 Jan 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20x), 100 mL	Not Kit Specific
Proinflammatory Panel 1 (human) Control 1	Not Kit Specific
Proinflammatory Panel 1 (human) Control 2	Not Kit Specific
Proinflammatory Panel 1 (human) Control 3	Not Kit Specific

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		24 SEP 2018